

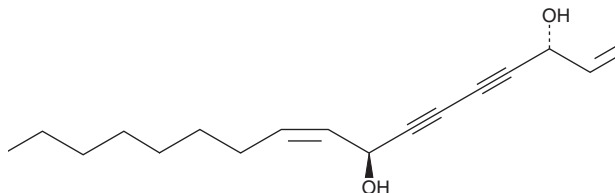
# PRODUCT INFORMATION



## (3R,8S)-Falcarindiol

Item No. 30919

CAS Registry No.: 225110-25-8  
Formal Name: (3R,8S,9Z)-1,9-heptadecadiene-4,6-diyne-3,8-diol  
Synonym: (+)-(3R,8S)-Falcarindiol  
MF:  $C_{17}H_{24}O_2$   
FW: 260.4  
Purity:  $\geq 98\%$   
Supplied as: A neat oil  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years  
Item Origin: Plant/*Notopterygium incisum*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

(3R,8S)-Falcarindiol is supplied as a neat oil. A stock solution may be made by dissolving the (3R,8S)-falcarindiol in the solvent of choice, which should be purged with an inert gas. (3R,8S)-Falcarindiol is soluble in ethanol, DMSO, and chloroform.

### Description

(3R,8S)-Falcarindiol is a polyacetylene that has been found in *N. incisum* and has diverse biological activities.<sup>1-3</sup> It is active against *M. tuberculosis* strain H37Ra (MIC = 6.25  $\mu\text{g/ml}$ ).<sup>1</sup> (3R,8S)-Falcarindiol is cytotoxic to MCF-7, H1299, and HepG2 cells ( $\text{IC}_{50}\text{s}$  = 29.4, 22.1, and 23.6  $\mu\text{M}$ , respectively).<sup>2</sup> It has algicidal activity against *H. circularisquama* when used at concentrations of 2 and 4  $\mu\text{M}$ .<sup>3</sup>

### References

1. O'Neill, T., Johnson, J.A., Webster, D., *et al.* The Canadian medicinal plant *Heracleum maximum* contains antimycobacterial diynes and furanocoumarins. *J. Ethnopharmacol.* **147**(1), 232-237 (2013).
2. Zheng, X., Zheng, X., Zhang, C., *et al.* Cytotoxic polyacetylenes isolated from the roots and rhizomes of *Notopterygium incisum*. *Chinese Chem. Lett.* **30**(2), 428-430 (2019).
3. Tamura, S., Ohno, T., Hattori, Y., *et al.* Establishment of absolute stereostructure of falcarindiol, algicidal principle against *Heterocapsa circularisquama* from *Notopterygium* Rhizoma. *Tetrahedron Lett.* **51**(11), 1523-1525 (2010).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/28/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM